
BD Accuri™ C6 Flow Cytometer

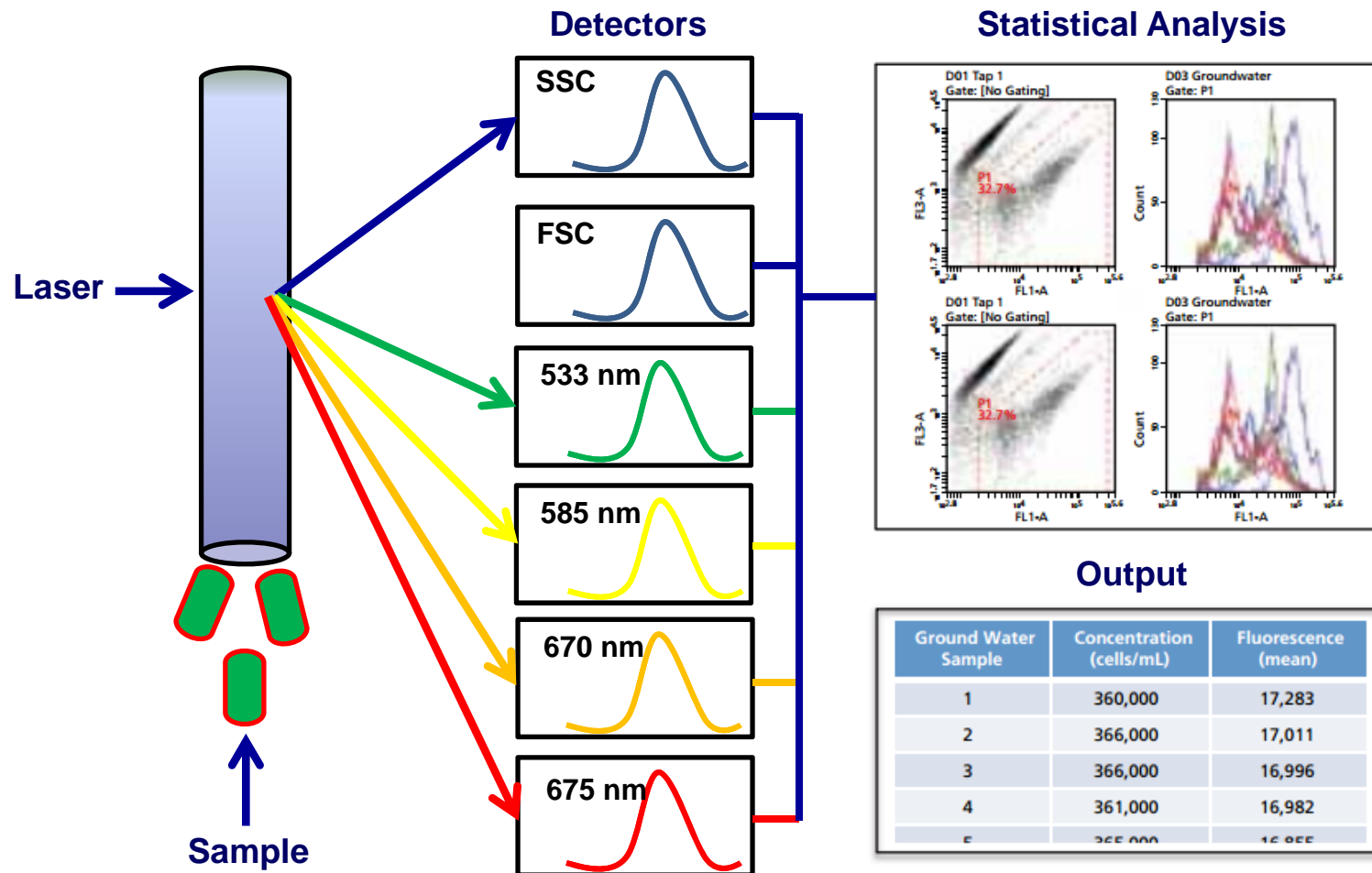
Mirko Corselli, PhD
BD Biosciences
Senior Scientist

The BD Accuri C6 Flow Cytometer System

- An affordable, full-featured, easy-to-use flow cytometer
- Two lasers and six detectors



BD Accuri C6 Flow Cytometer



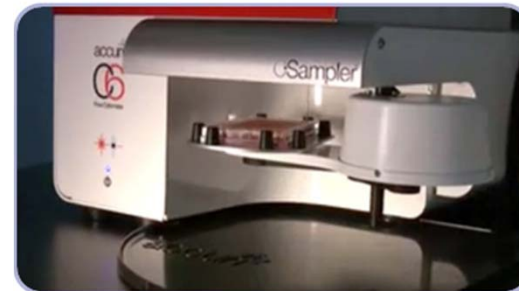
Advantages of Pre-optimized Detector Settings



- **Greatly reduces the risk of lost data due to improper setup**
- **Saves time and sample**
- **No specialist training or dedicated operator required**
- **Predictable, reproducible analysis relative to the sample type and application**

Enhanced Sample Handling

- **Direct volume measurement**
- **Many types of sample tubes may be used.**
 - Flow cytometry tubes
 - Microcentrifuge tubes
 - Ninety-six-well plates with the BD CSampler™ accessory
- **Open system conducive to kinetic studies**
- **BD CSampler™ accessory for automated sample introduction**



Intuitive Software



Sample Grid

Cytometer Status

Fluidics Controls

Run Criteria

Real-Time Updates

The screenshot displays the BD Accuri C6 software interface, divided into several functional areas:

- Collect Tab:** Shows a sample grid for 'A4' with parameters 'HPB CD3-F CD4, CD45, CD8'. The grid has columns 1-12 and rows A-H. Cell A4 is highlighted.
- Analyze Tab:** Contains six flow cytometry plots (Plot 1 to Plot 6) showing various parameters like SSC-A vs FSC-A, CD4 PE-A vs CD3 FITC-A, CD45 PE-CY7-A vs CD3 FITC-A, CD4 PE-A vs CD4 APC-A, and CD8 APC-A vs CD3 FITC-A. Each plot includes a gate (P1, Q1-Q5) and percentage values.
- Statistics Tab:** Displays two tables of data for the analyzed plots.
- Control Panels:** Includes 'Run Limits' (125,000 events), 'Fluidics' (Flow Rate 14 µL/min, Core Size 10 µm), and 'Threshold' (80,000 on FSC-H).
- Buttons:** 'ADD to A4', 'Set Color Compensation', 'Backflush', 'Unglog', and 'Delete Sample Data'.

	Count	Volume (µL)	% of This Plot	% of All	Mean CD3 FITC-A	Mean CD4 PE-A	CV CD3 FITC-A	CV CD4 PE-A
This Plot	99,183	0.0	100.0%	43.3%	23,875.3	1,762.1	62.0%	62.0%
Q3-UL	147	0.0	0.1%	0.1%	960.4	1,607.0	53.3%	53.3%
Q3-UR	59,194	0.0	59.7%	25.8%	32,131.1	2,933.7	28.5%	28.5%
Q3-LR	18,387	0.0	18.5%	8.0%	24,701.0	22.5	29.2%	29.2%

	Count	Volume (µL)	% of This Plot	% of All	Mean CD3 FITC-A	Mean CD45 PE-CY7-A	CV CD3 FITC-A	CV CD45 PE-CY7-A
This Plot	102,719	0.0	100.0%	44.8%	23,071.3	38,996.7	65.7%	65.7%
R1	99,183	0.0	96.6%	43.3%	23,875.3	40,270.2	62.0%	62.0%
Q1-UL	21,694	0.0	21.1%	9.5%	540.5	32,690.2	45.5%	45.5%
Q1-UR	77,625	0.0	75.6%	33.9%	30,356.9	42,417.0	30.7%	30.7%
Q1-LR	3,380	0.0	3.3%	1.5%	447.0	1,141.0	76.5%	76.5%
Q1-LL	20	0.0	0.0%	0.0%	8,449.7	2,092.6	97.0%	97.0%

Histogram, Dot Plot, and Density Plot Display Area

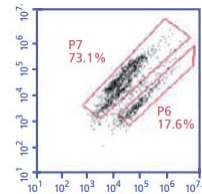
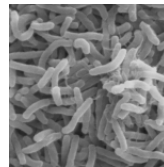
Analysis and Gating Tools

Plot Statistics

A Versatile Instrument for Broad Applications

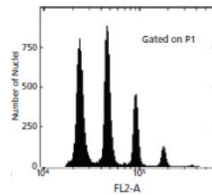
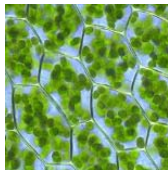


Microbiology



- Aquatic microbiome analysis
- Biofuel research
- Bacteria viability and concentration

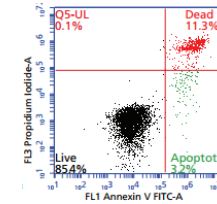
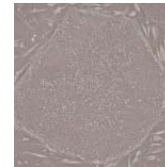
Plant Biology



- DNA content

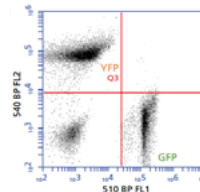
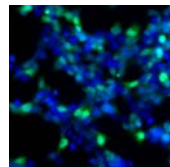


Cell Biology



- Apoptosis
- Proliferation
- Immunophenotyping

Fluorescent Protein Analysis



- GFP, YFP
- mCherry®, RFP
- mOrange®, dTomato®

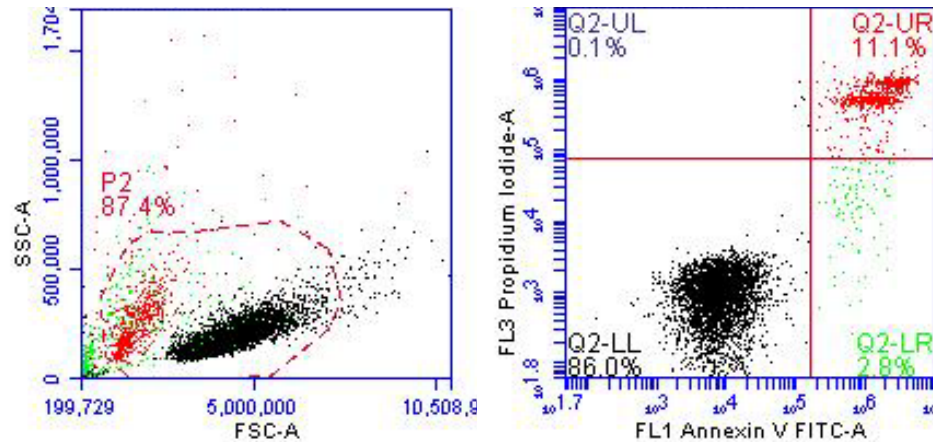
Flow Cytometry within Reach™

The BD Accuri™ C6 Personal Flow Cytometry Tour

Apoptosis: Annexin V



DMSO



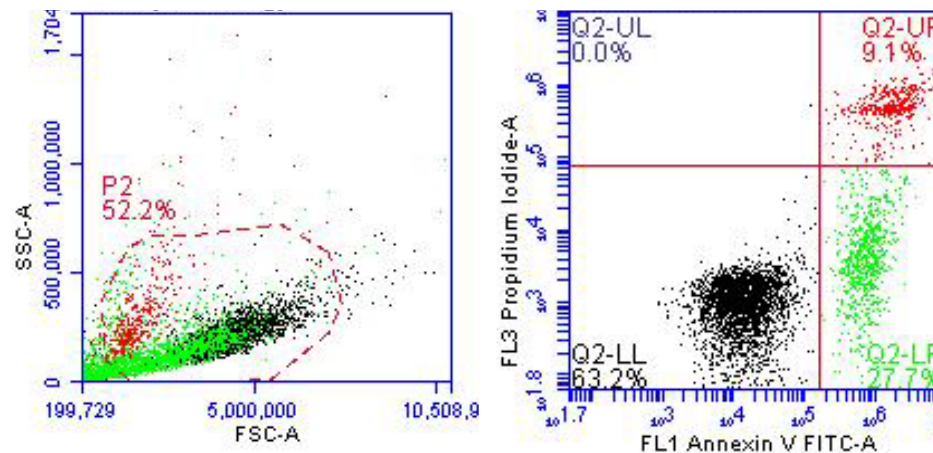
Live cells:

**Annexin V Negative
PI Negative**

Early apoptotic:

**Annexin V Positive
PI Negative**

Camptothecin
6 μ M

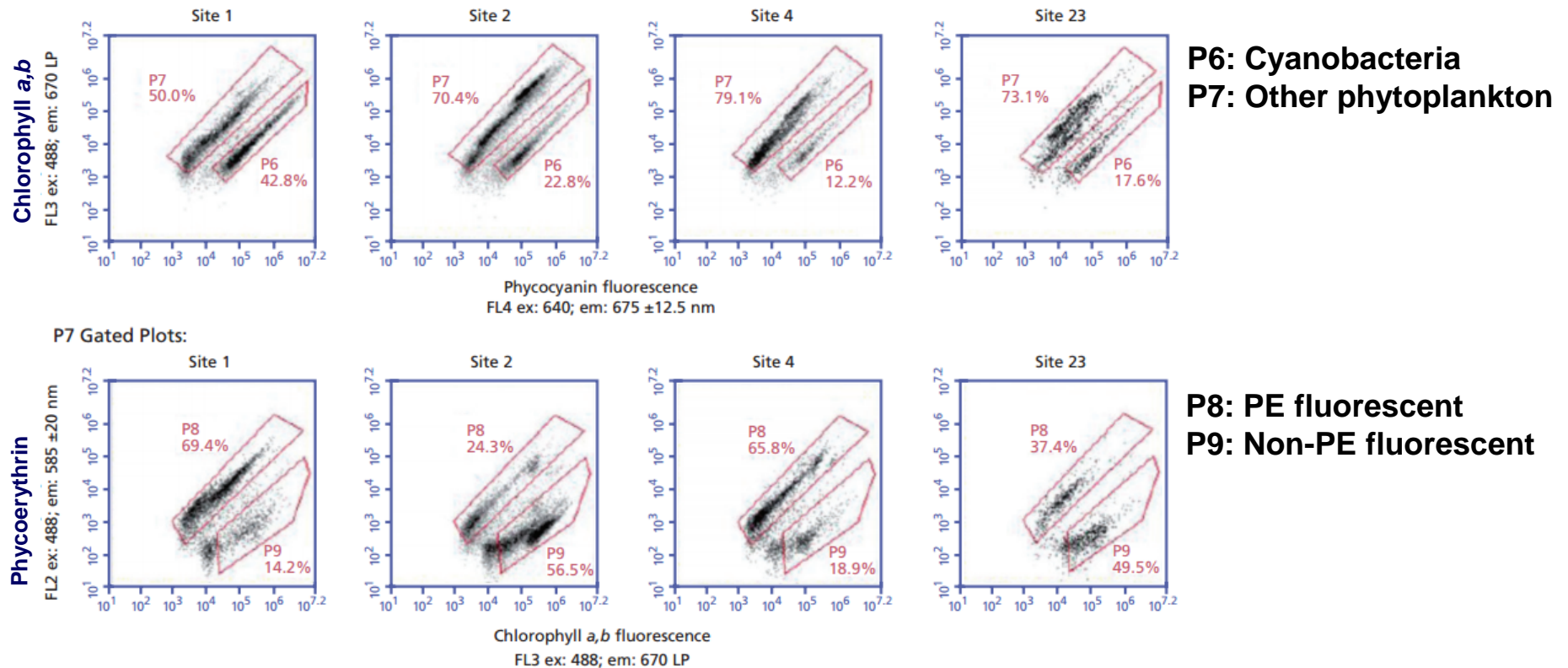


Late apoptosis/dead:

**Annexin V Positive
PI Positive**

Jurkat cells were treated with DMSO or camptothecin for 4 hours.

Analysis of Aquatic Samples from Saginaw Bay: Fluorescence



Four types of phytoplankton were identified by fluorescence characteristics.

Data courtesy of J.D. Bressie, PhD, NOAA, Seattle, WA

Kits and Templates on the BD Accuri C6



Category	Product Information Sheet	Brand	Kit	Cat. No.	Template
Cell Biology	BD Apoptosis Kits and Templates	BD Pharmingen™	Annexin V FITC Apoptosis Detection Kit II	556570	Download
		BD Pharmingen™	Annexin V PE Apoptosis Detection Kit I	559763	Download
		BD™	MitoScreen (JC-1) Kit	551302	Download
		BD Pharmingen™	Caspase-3 PE Assay Kit	550914	Download
		BD Pharmingen™	Caspase-3 FITC Assay Kit	550480	Download
	BD Cell Cycle and DNA Kits and Templates	BD Cycletest™ Plus	DNA Reagent Kit	340242	Download
		BD Pharmingen™	FITC BrdU Flow Kit	559619	Download
		BD Pharmingen™	APC BrdU Flow Kit	552598	Download

BD Accuri C6 Promotion and Personal Flow Cytometry Tour



**BEST
IN CLASS**
meets
BEST
TIME TO BUY

10% off
BD Accuri™ C6

40% off
BD Reagents

The graphic features a dark blue background with two overlapping circles. The left circle contains a white BD Accuri C6 flow cytometer with a red top and is labeled '10% off' and 'BD Accuri™ C6'. The right circle contains various reagent bottles and boxes and is labeled '40% off' and 'BD Reagents'. A white plus sign is positioned between the two circles.

Note:

US Region Only

Promotion Period: Oct 1, 2014 – Dec 31, 2014

The BD Accuri C6 Personal Flow Cytometry Tour

- Introduction to Flow Cytometry
- Cancer and Cell Biology Applications
- Microbial Analysis

Flow Cytometry within Reach™

The BD Accuri™ C6 Personal Flow Cytometry Tour

For Additional Information...



BD Biosciences

Sign In



Home / Instruments and Software / BD Accuri C6

BD ACCURI C6 [Overview](#) [Features](#) [Applications](#) [Products](#) [Sample Data](#) [Resources & Tools](#)

Supports cell analysis for up to six parameters

The BD Accuri™ C6 makes the analytical power of flow cytometry more accessible with ease-of-use and affordability. Its compact footprint and portable weight make it a valuable personal use tool for both novice and experienced researchers who want a cytometer to be easily available when and where they need it.

The system features an intuitive software interface, software templates, and reagent kits that guide users new to flow cytometry through workflows for popular applications.



[REQUEST DEMO](#)

[GET QUOTE](#)



RESEARCH APPLICATIONS

Simplify setup and analysis for immunophenotyping, apoptosis, cell cycle, microbial counting, and intracellular cytokines.



SAMPLE DATA

View experiment results from a wide range of applications including gene expression, cell cycle, and DNA analysis.

[CONTACT US](#)

Key Resources

[Brochure](#)
[Filter Guide](#)
[Technical Specifications](#)
[Kits and Templates Catalog](#)
[All Resources »](#)

Community

[BD Accuri News](#)

Related Links

[Services](#)

Related Products

[Multicolor Reagents](#)
[Bead-Based Immunoassays](#)
[Apoptosis, Cell Cycle, and Cell Proliferation](#)

More Information

[Ask BD](#)
[News and Events](#)
[Sign Up for Email Updates](#)

www.bdbiosciences.com/resources/accuri

Technical Support:

Ph: 877-232-8995, Prompt 3, 2

email: ResearchApplications@bd.com

Class 1 Laser Product.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

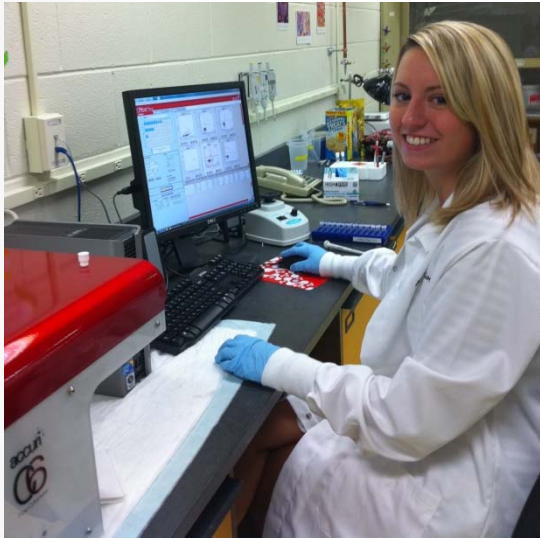
Living Colors® (including mCherry, mOrange, and dTomato dyes) is a registered trademark of Clontech.

BD, BD Logo and all other trademarks are property of Becton, Dickinson and Company. © 2014 BD

23-16980-00



Flow cytometry in undergraduate education



Whitney Edwards



**Samantha Scott, Talbot Weston, Sarah Murphy,
Melanie Gubbels Bupp – Autumn Immunology
Conference 2014**

Melanie Gubbels Bupp, Assistant
Professor of Biology

General Advice on using cytometry in higher education

- Use the technology for projects that allow students to...
 - more firmly grasp basic concepts
 - test something of interest to them
- Only tell them what they need to know about the technology WHEN they need to know it.
 - Front-loading a lot of technical information is overwhelming and often, ineffective



Examples of using cytometry at R-MC

- **Peritonitis Lab** in Immunology course for biology majors [simplified version taught in a freshman mixed (majors/non-majors) lab]
- **Phagocytosis Lab** in Immunology course for biology majors
- **Independent research projects** for undergraduate students



Introduce students to the lab

Thioglycollate-Induced Peritonitis: Recruitment of leukocytes from the circulation, and their subsequent influx into the sites of inflammation is critical for host defense and wound healing. This is a multistep process, which is regulated, in part, by adhesion molecules and chemokines that are upregulated during inflammation. An intra-peritoneal injection of thioglycollate generates local inflammation and initiates the migration of inflammatory cells to the site of inflammation. Thus, thioglycollate-induced peritonitis in mice mimics an acute inflammatory response in the peritoneum.

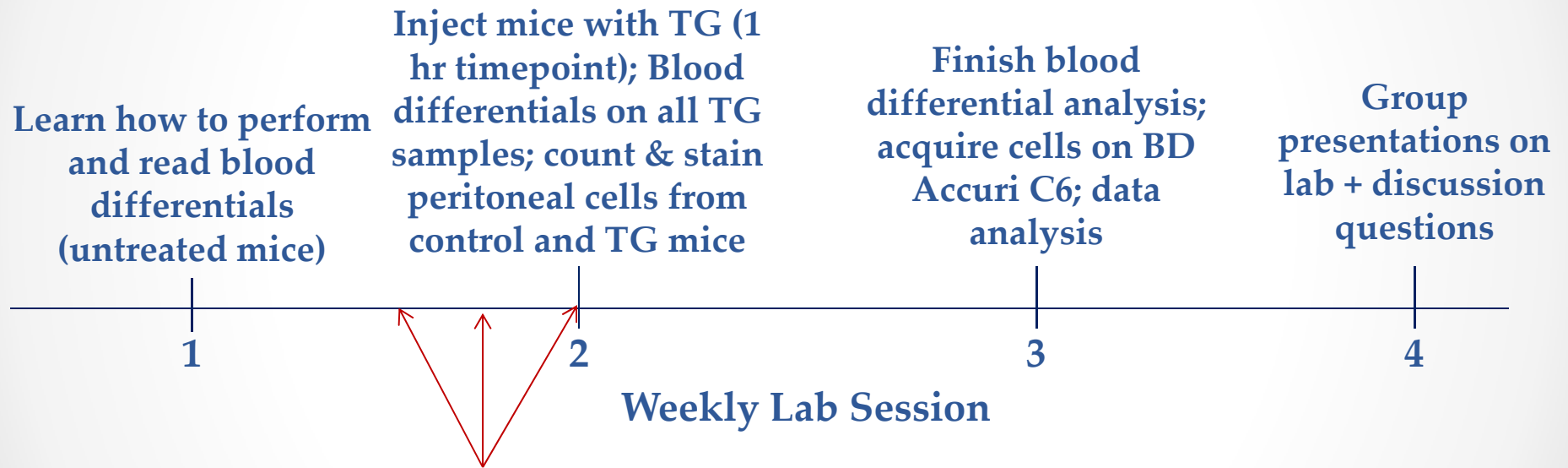


Learning Objectives

- Be able to identify particular white blood cells (lymphocytes, monocytes, neutrophils) in a blood smear.
- Explain how and why blood differentials are taken
- Describe the chronological order in which immune cells arrive at sites of inflammation, such as the peritoneal cavity in thioglycollate-treated mice.



Peritonitis Lab Timeline



Inject mice with 6% thioglycollate (TG) in PBS or control PBS to induce peritoneal inflammation

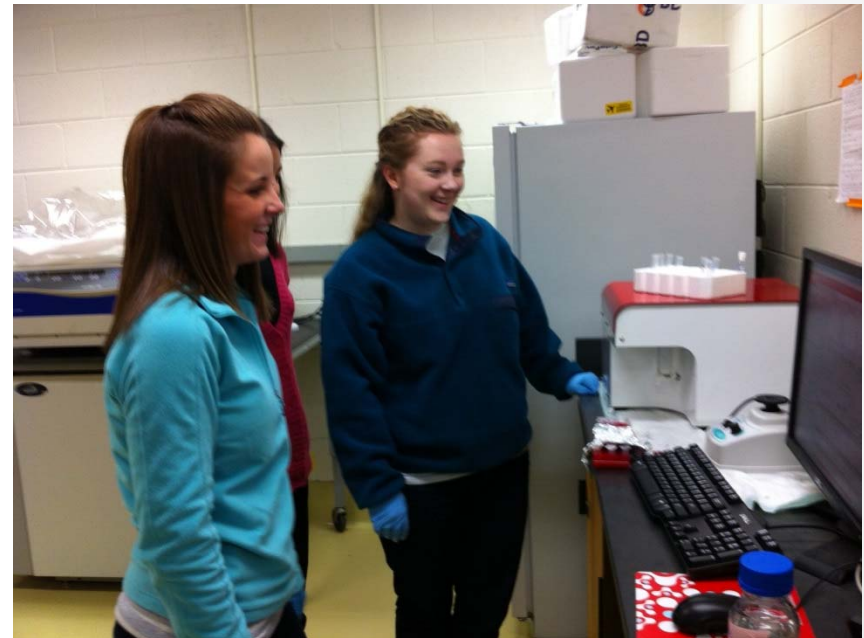


R-MC Immunology students acquiring their samples

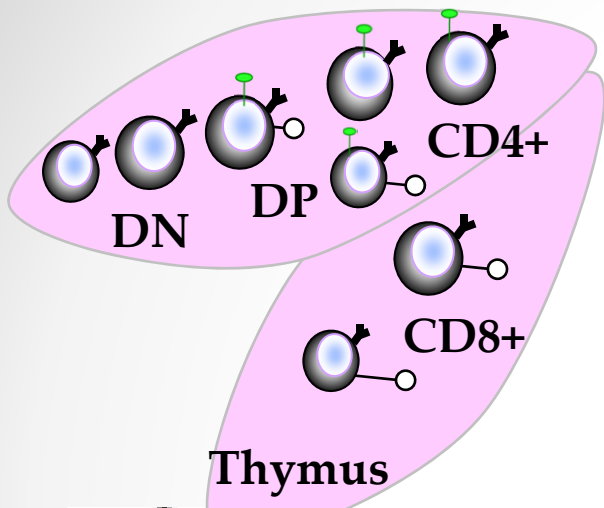
Victoria Robinson



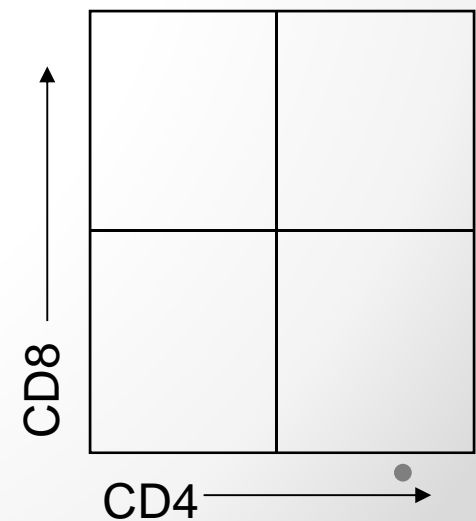
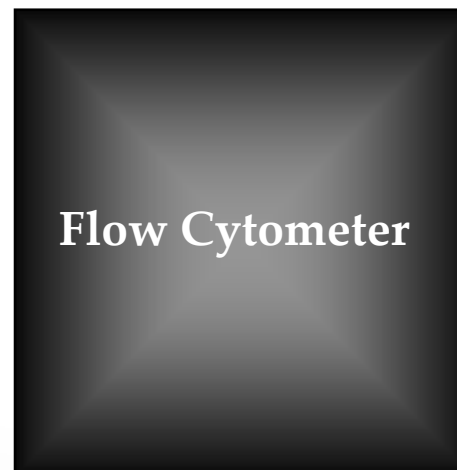
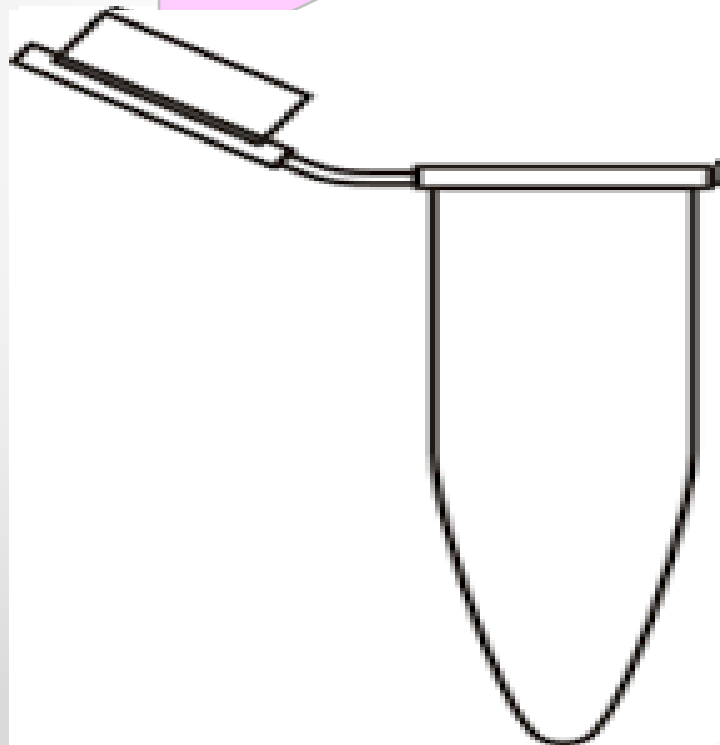
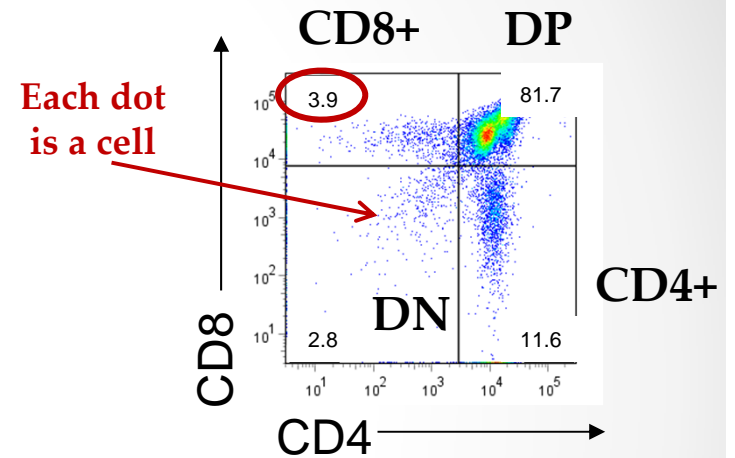
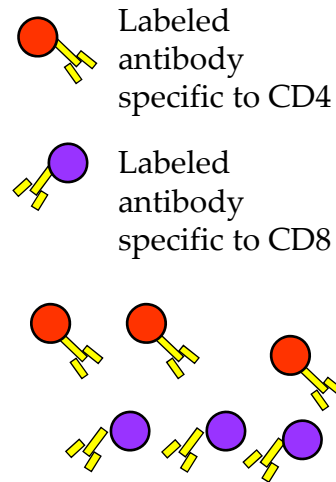
**Lauren Philips, Jane Oh, &
Casey Kaufman**



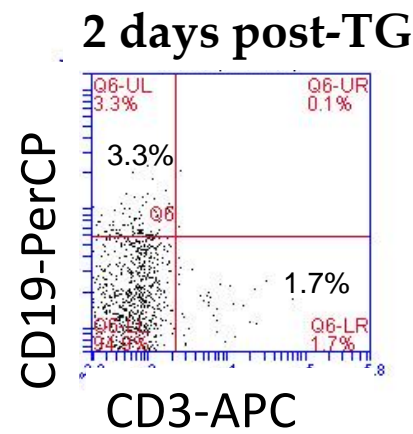
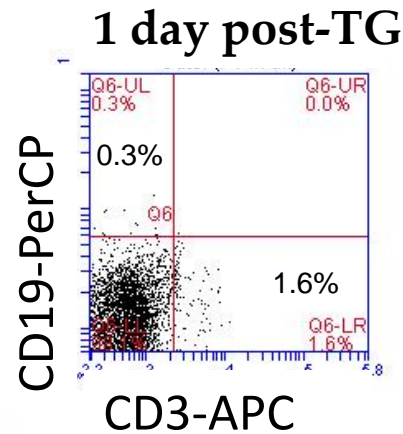
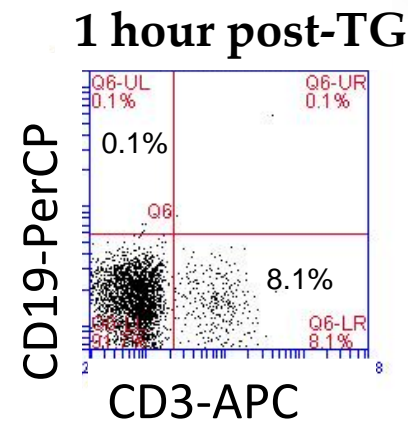
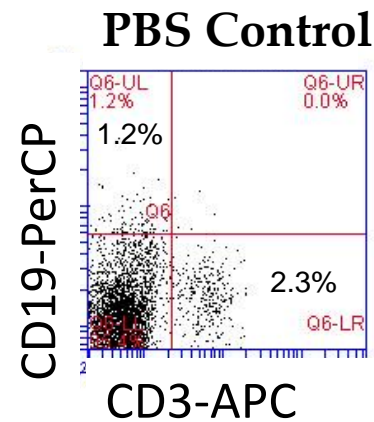
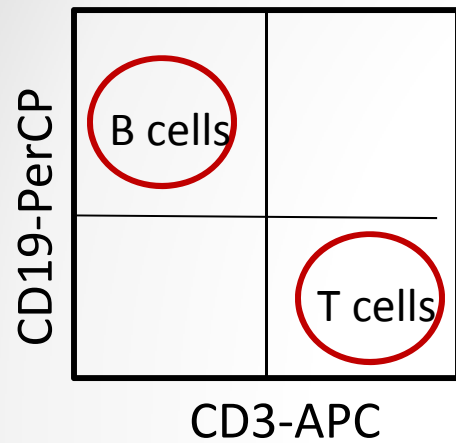
Explaining how cytometry works



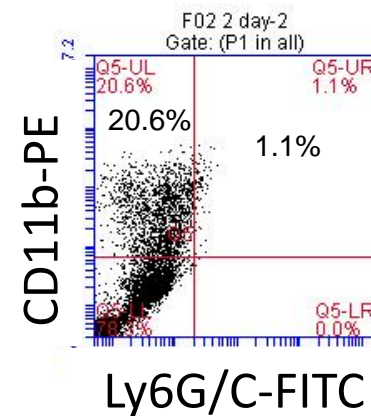
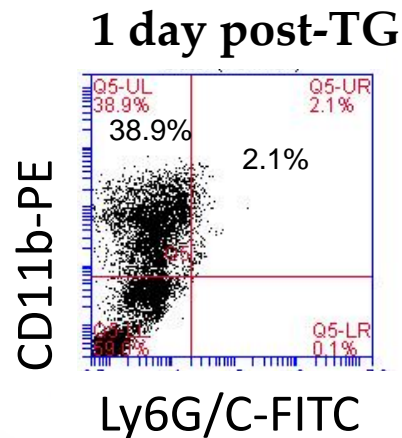
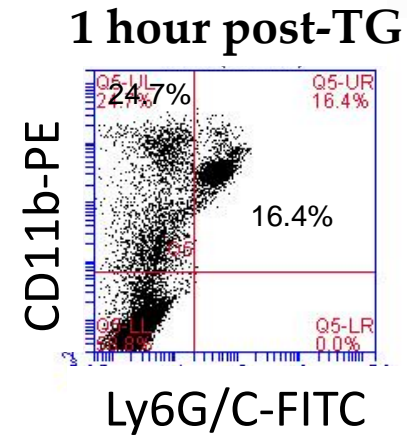
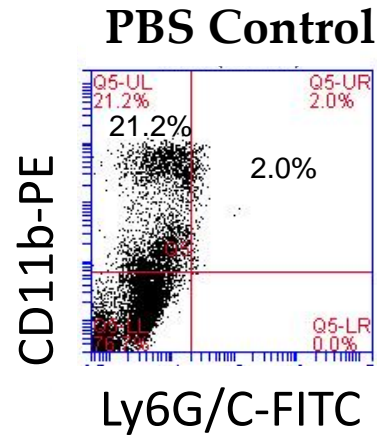
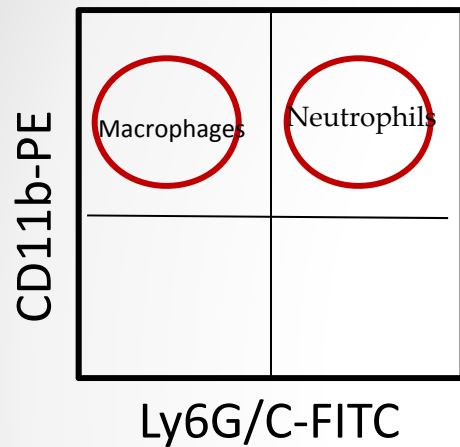
works



Example student data



Example student data




Example discussion questions for student peritonitis presentations

- What types of junctions exist between endothelial cells lining blood vessel walls in the non-inflamed, resting condition? How do these junctions change during local inflammation?
- Relate your findings from the blood differentials with your findings from the flow cytometry data. Do the two sets of data “paint the same picture”? Why or why not?
- Map the route newly developed neutrophils must take to enter the inflamed peritoneal cavity. Begin in the bone marrow and end with the peritoneal cavity.

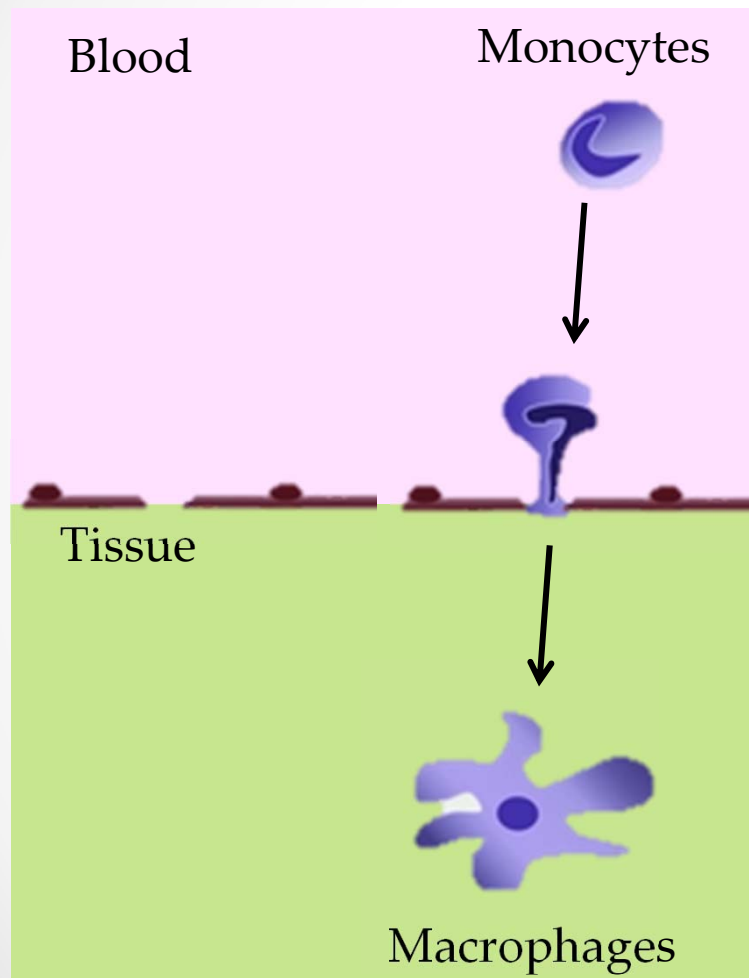


Examples of using cytometry at R-MC

- **Peritonitis Lab** in Immunology course for biology majors [simplified version taught in a freshman mixed (majors/non-majors) lab]
-  **Phagocytosis Lab** in Immunology course for biology majors
- **Independent research projects** for undergraduate students



Introduce students to the lab



Learning Objectives

- Compare and contrast monocytes and macrophages
- Evaluate flow cytometry data regarding the ability of cells to phagocytose fluorescently tagged antigens
- Be able to design an experiment to test the impact of various compounds on phagocytosis



Phagocytosis Lab Timeline

Learn how to culture THP-1 cells; design experiment to test effect of particular substance on phagocytosis (*in vitro*)

Incubate THP-1 cells with FITC-latex beads, and acquire samples on BD Accuri C6; data analysis

Group presentations on lab + discussion questions

1

2

3

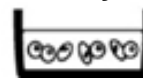
Weekly Lab Session

Set up THP-1 cultures in 24 well plates +/- PMA and +/- test substance

THP-1 monocytes

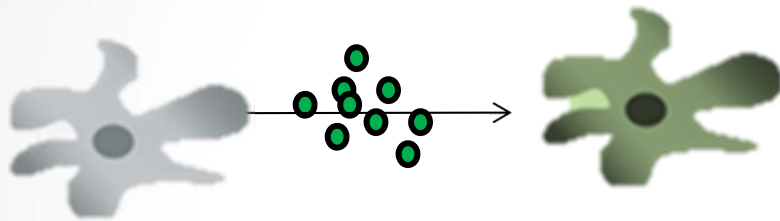
+PMA
24 h

THP-1 macrophages

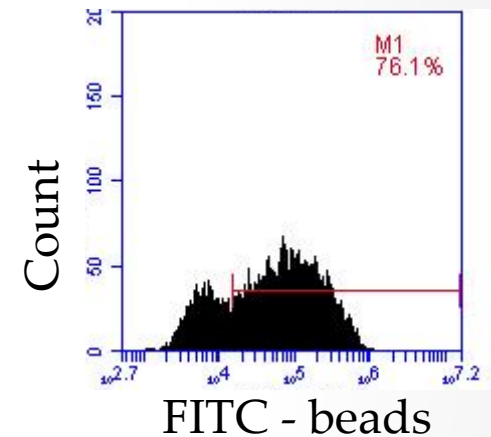
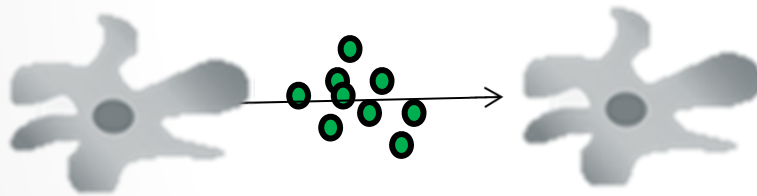


Explaining how cytometry works -- histograms

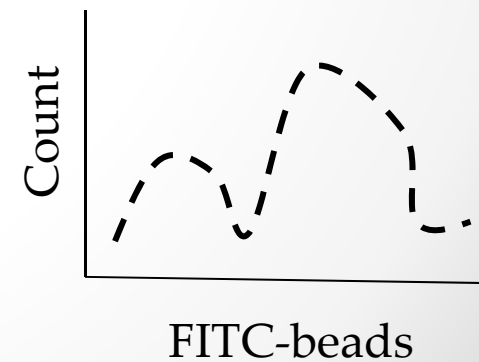
Lots of phagocytosis



Very little phagocytosis



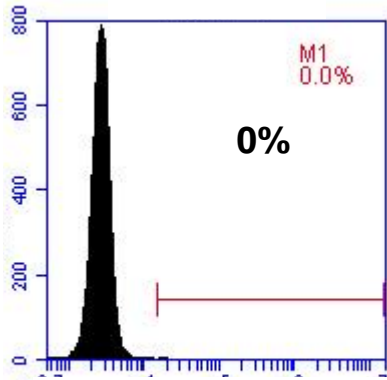
Flow Cytometer



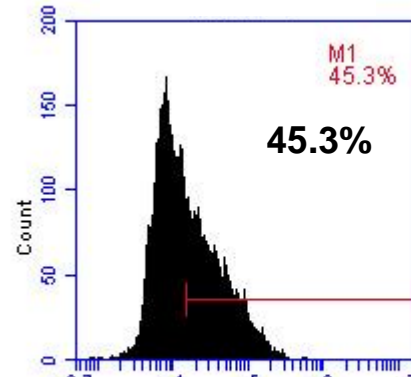
Example student data

0 PMA

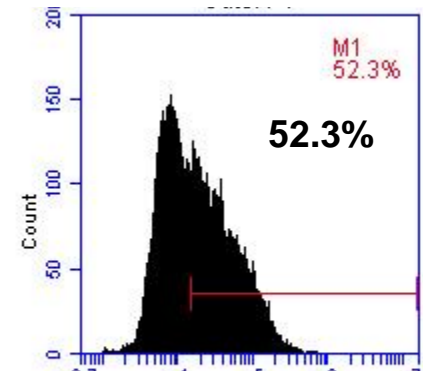
THP cells



THP cells +
FITC-beads

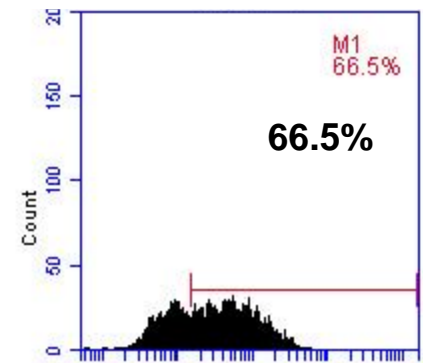
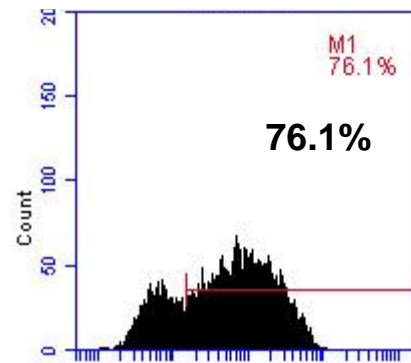
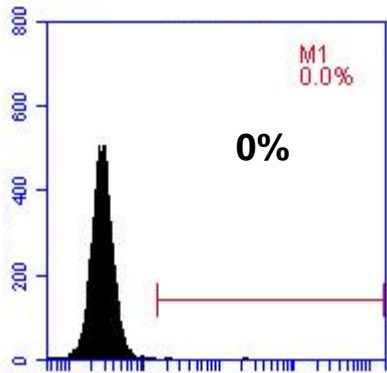


THP cells +
FITC-beads +
Vitamin C



+ PMA

Count



Examples of using cytometry at R-MC

- **Peritonitis Lab** in Immunology course for biology majors [simplified version taught in a freshman mixed (majors/non-majors) lab]
- **Phagocytosis Lab** in Immunology course for biology majors
- **Independent research projects** for undergraduate students

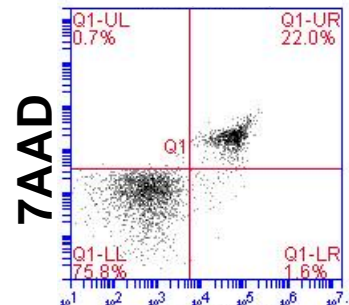


Students at R-MC also use the cytometer in independent research projects

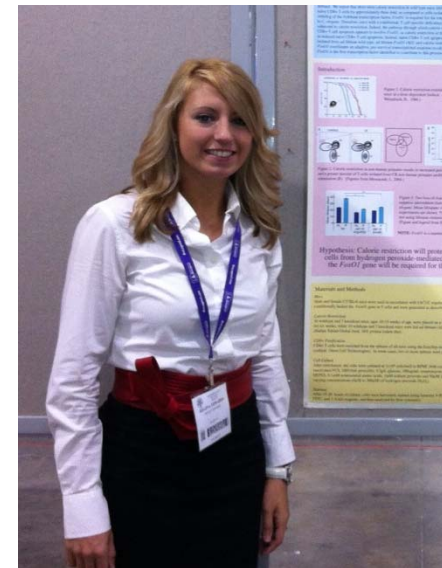
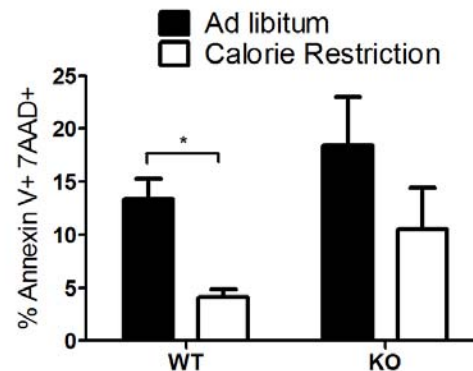


Seth Litvin

Apoptosis Markers



Annexin V - PE



Whitney Edwards



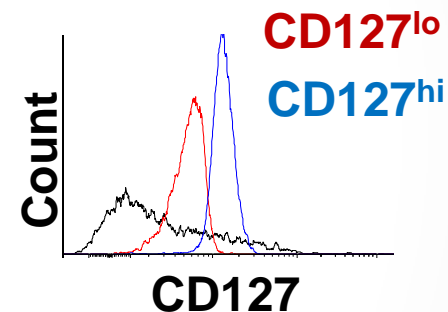
Students at R-MC also use the cytometer in independent research projects

Sarah Murphy



Isolate Naïve CD8+ T cells from Donor mice

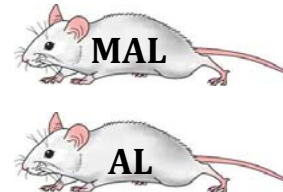
Sort them on CD127 expression



Adoptively Transfer cells into recipients

**CFSE^{hi}
CD127^{lo}
2.5 x 10⁶**

**CFSE^{lo}
CD127^{hi}
2.5 x 10⁶**

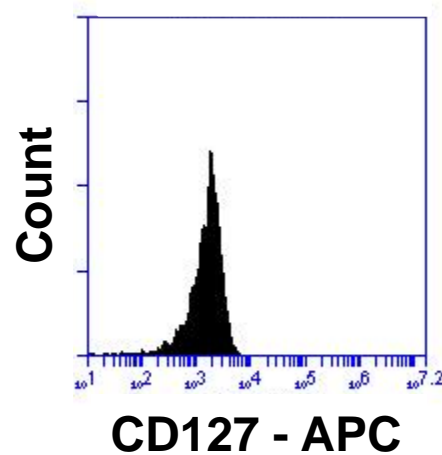
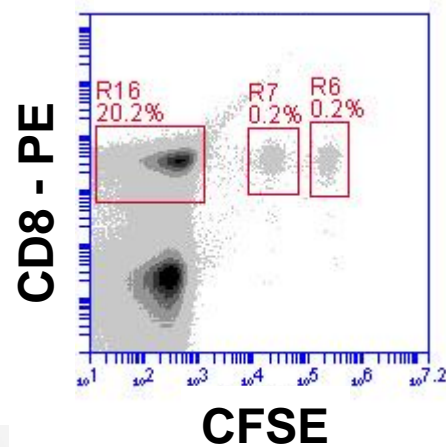


1 week

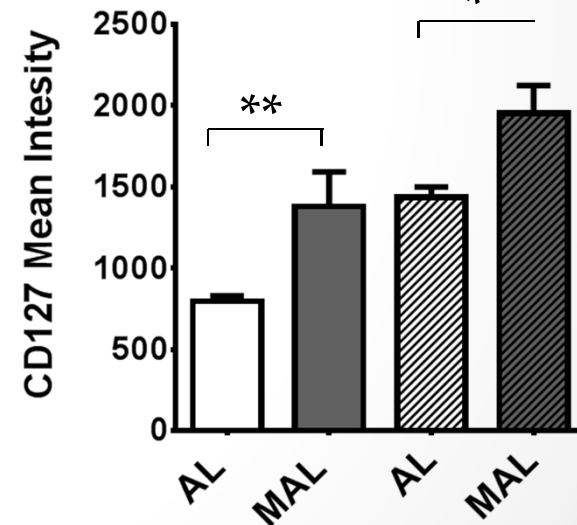
Assess total numbers and CD127 expression of CFSE+ cells by flow cytometry



Students at R-MC also use the cytometer in independent research projects



Naïve CD8⁺ CFSE⁺ T Cells
in recipient LNs



Donor cells: CD127^{low} CD127^{high}



Thank you to R-MC students participating in labs and independent research

- Whitney Edwards
- Erica Horseman
- Rebecca Davis
- Seth Litvin
- Victoria Robinson
- Josh Anoff
- Brittany Mihalcoe
- Alex Koppleman
- Samantha Scott
- Sarah Murphy
- Talbot Weston



**Talbot Weston, Sarah Murphy,
and Samantha Scott**



Integrating *Microbial* Flow Cytometry Into Education

Tim W Overton

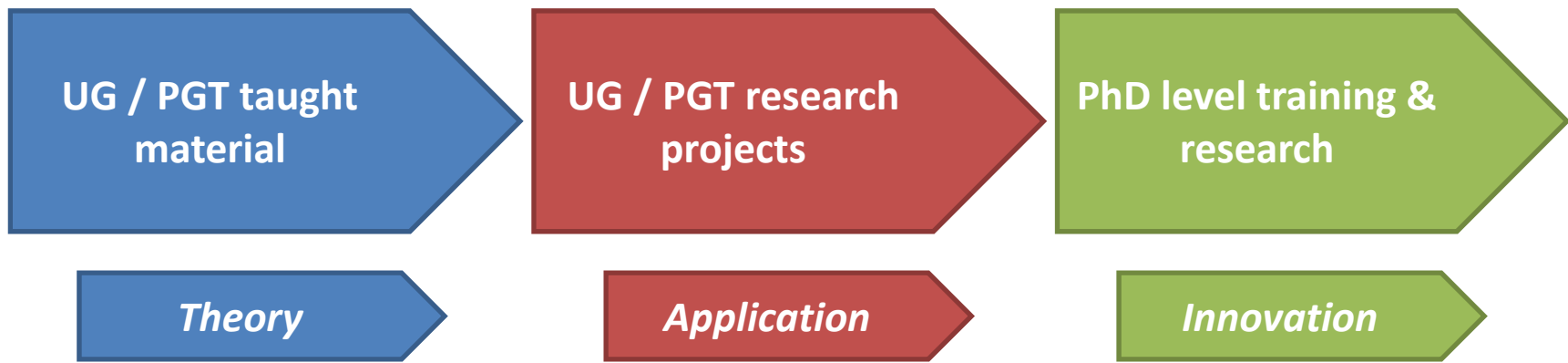
Bioengineering, School of Chemical Engineering

University of Birmingham

t.w.overton@bham.ac.uk @overtonlab

How can we get FCM applied in microbiology / microbial biotechnology?

- Collaboration between FCM specialists and microbiology / bioprocessing researchers
- Training at University level



Teaching fermentation

- MSc Biochemical Engineering

- Fermentation and cell culture for production of biopharmaceuticals

- Downstream processing of biopharmaceuticals

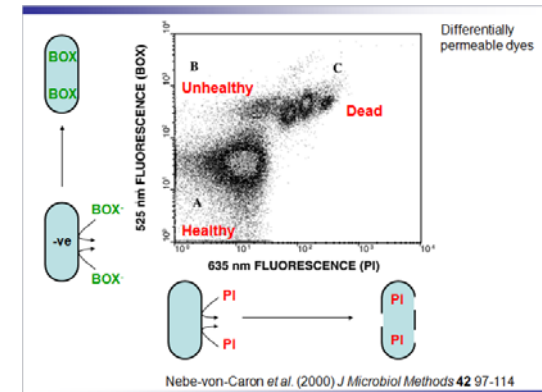
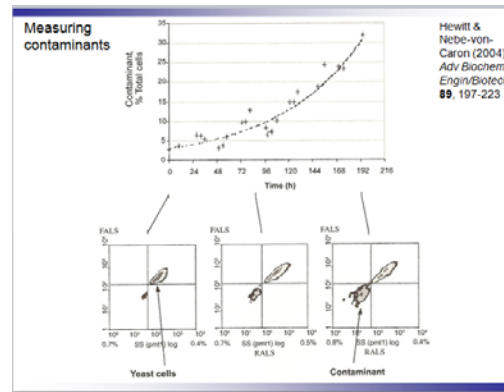
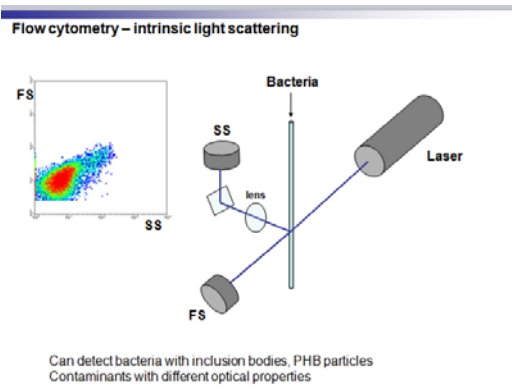
- Systems and synthetic biology

- Pharmaceutical, food and business themes

- Research project

Teaching analysis of fermentation

- Theory sessions on analytical techniques
 - Online versus offline
 - Real time / non-real time
 - Bulk versus single cell
 - Direct observations
- Theory and advantages of FCM

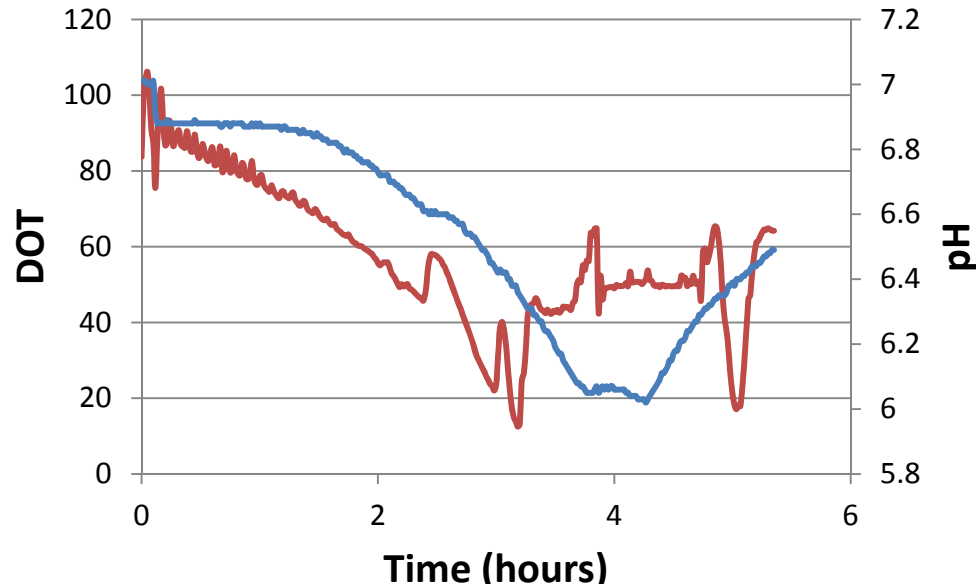


Lab-scale fermentation

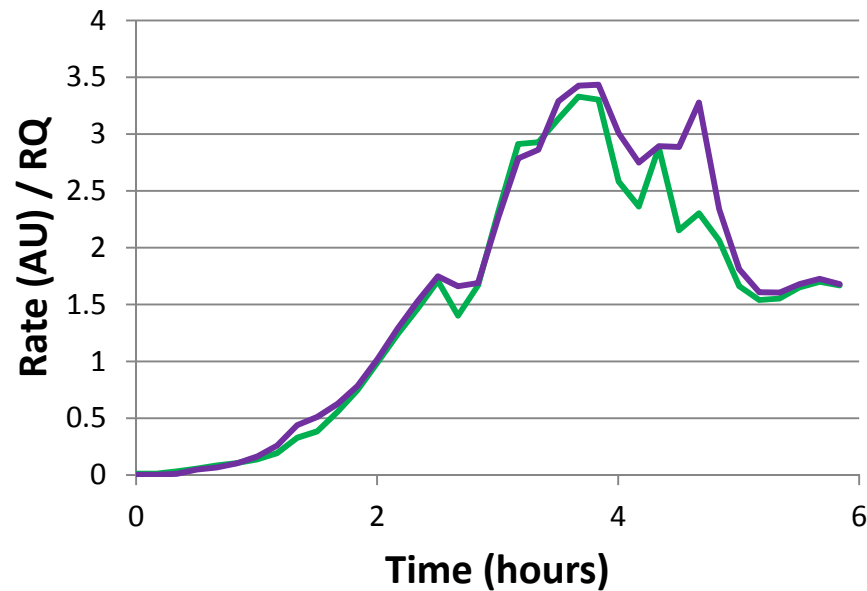
- Growth of *E. coli* in 5 litre bioreactors
- Analysis:
 - Traditional techniques: pH, DOT, offgas
 - Biomass measurements:
 - Optical density
 - Colony forming units
 - Dry cell weight
 - FCM



Fermentation data



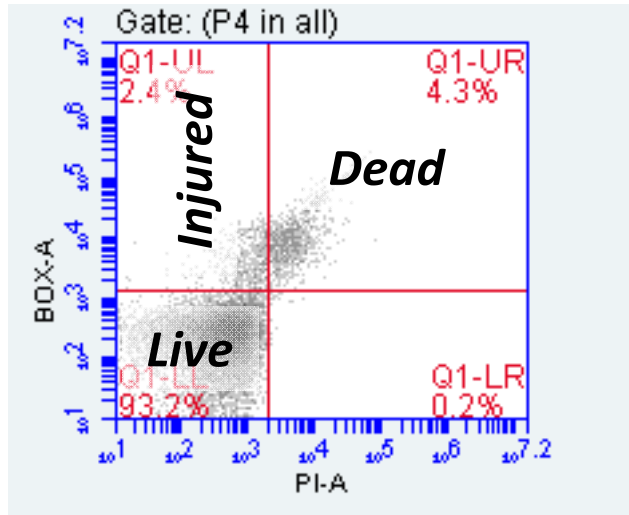
Traditional techniques:
Dissolved O₂ and **pH**
Basics of growth and physiology



O₂ consumption,
CO₂ evolution
Basics of growth and physiology

FCM with physiology dyes

1 hour



DiBac₄(3) (Bisoxanol; BOX)
Only enters depolarised cells
Stains depolarised cells green

Propidium Iodide (PI)

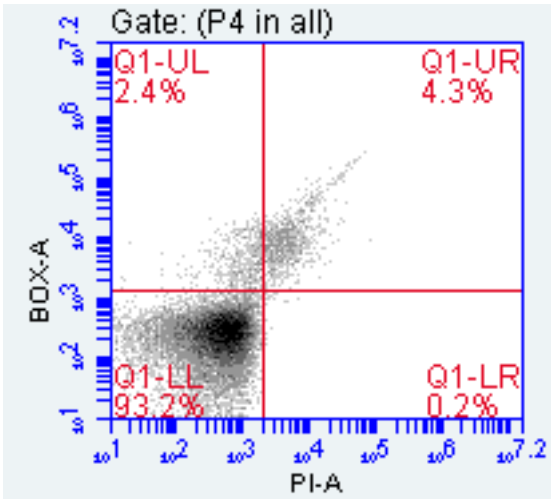
Enters cells through holes in wall
Stains dead cells red

Advantages:

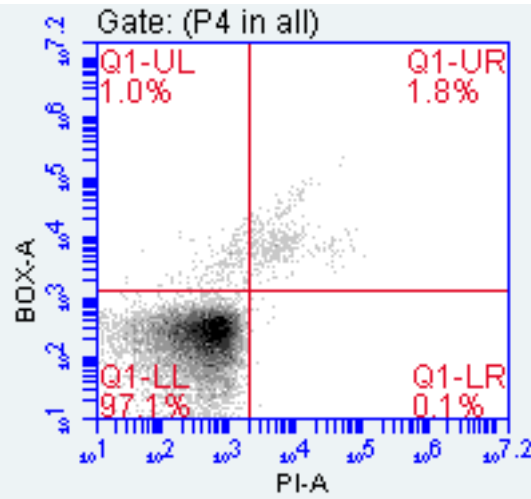
- Rapid method
- Allows monitoring of VBNC bacteria
- Allows counting of cells

FCM with physiology dyes

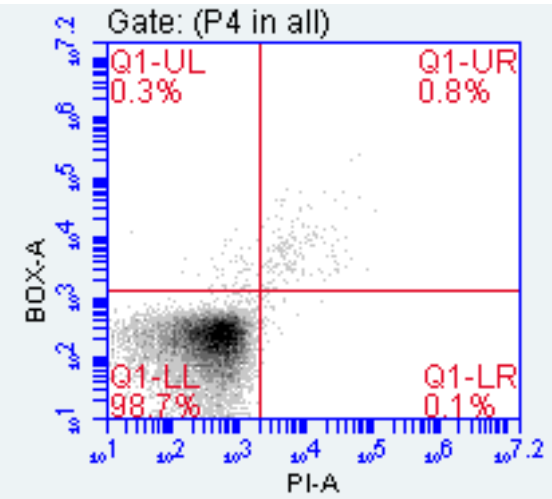
1 hour



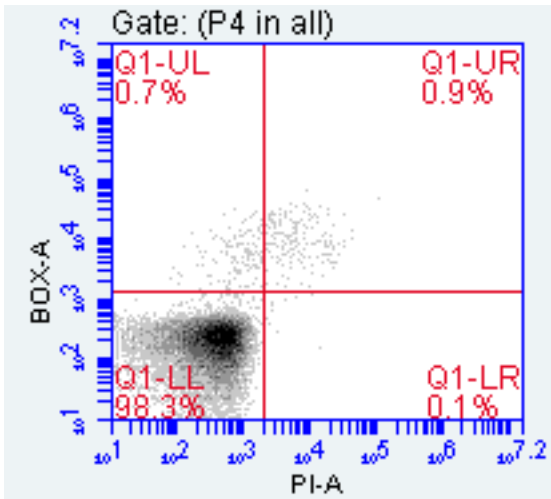
2 hours



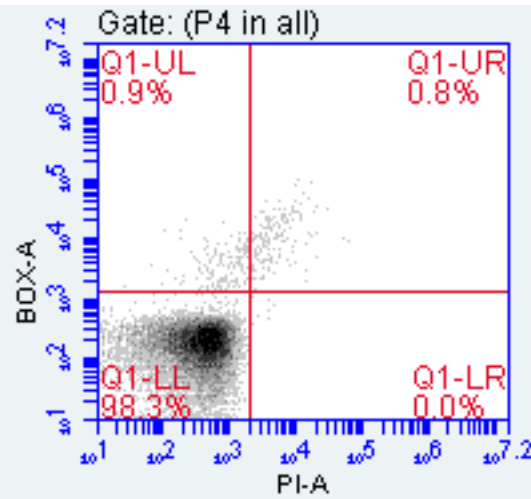
3 hours



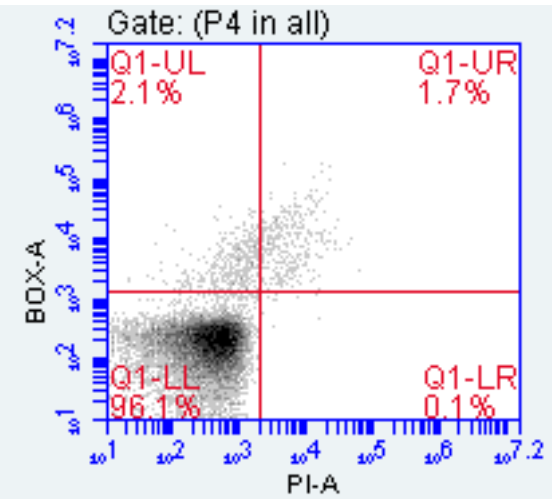
4 hours



5 hours



6 hours

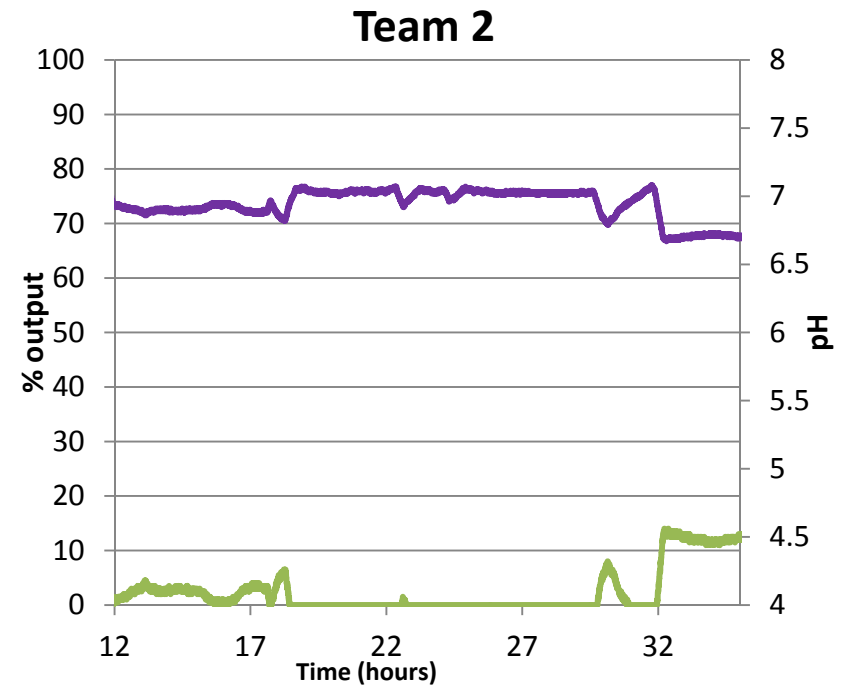
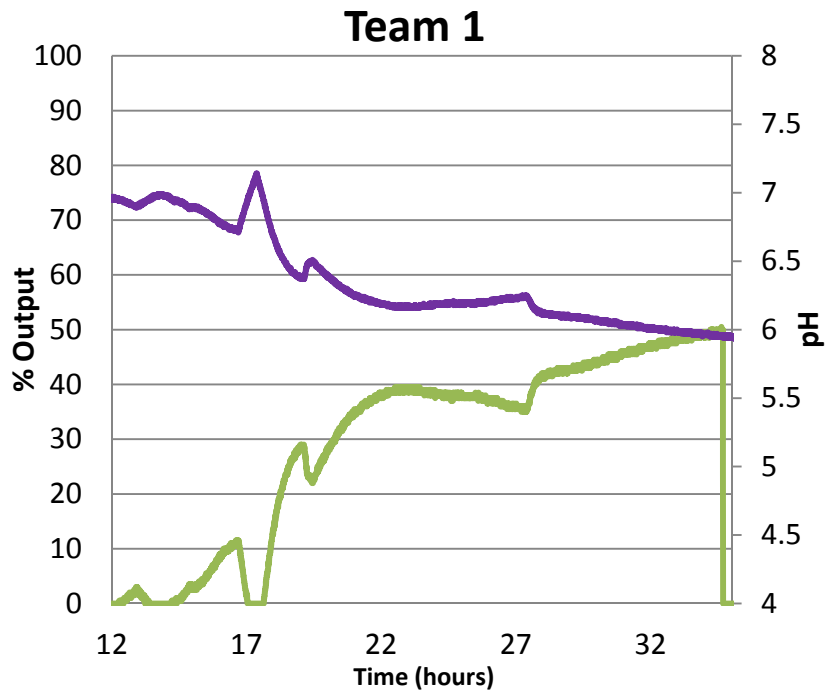


Pilot plant

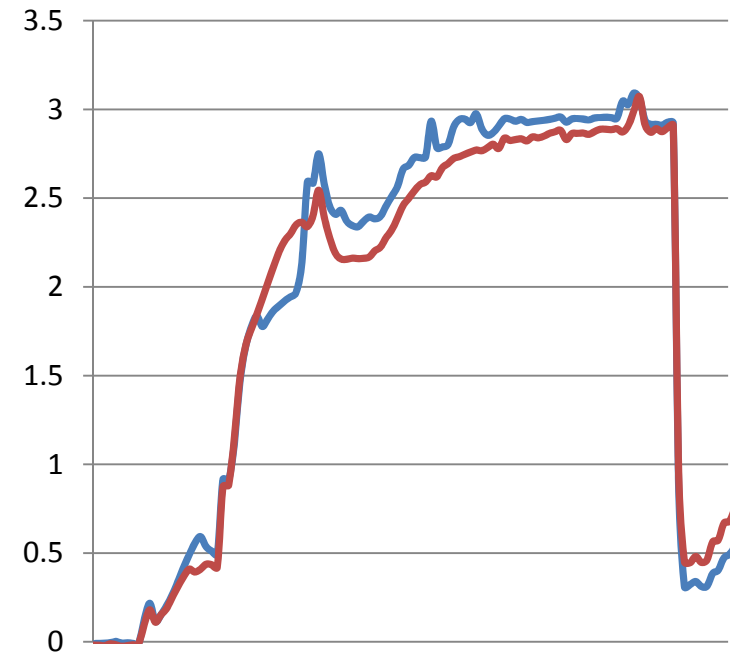
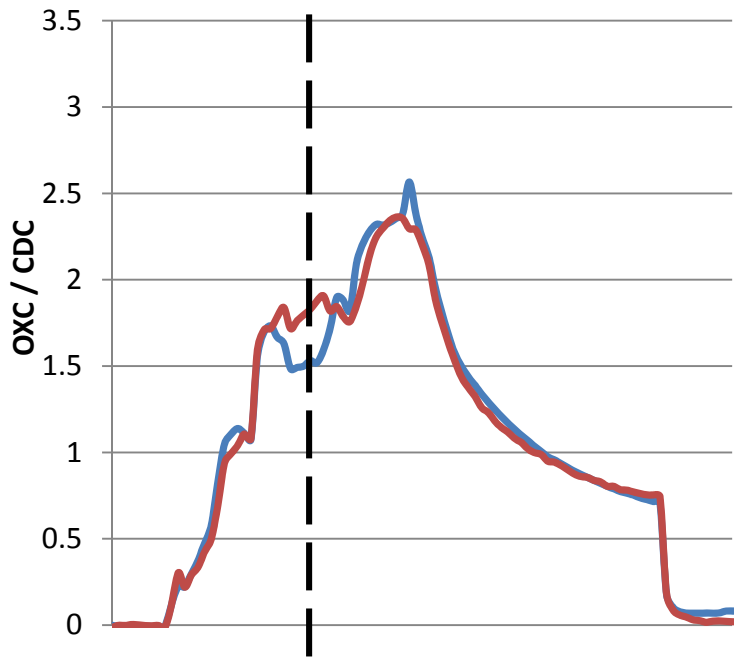
- Growth of *E. coli* in 150 litre vessels over the course of a week
- Fed-batch with glucose
 - Glucose feeding rate is critical to success
- Students encouraged to develop their own analysis methods and strategies – including FCM



Base addition rate & pH

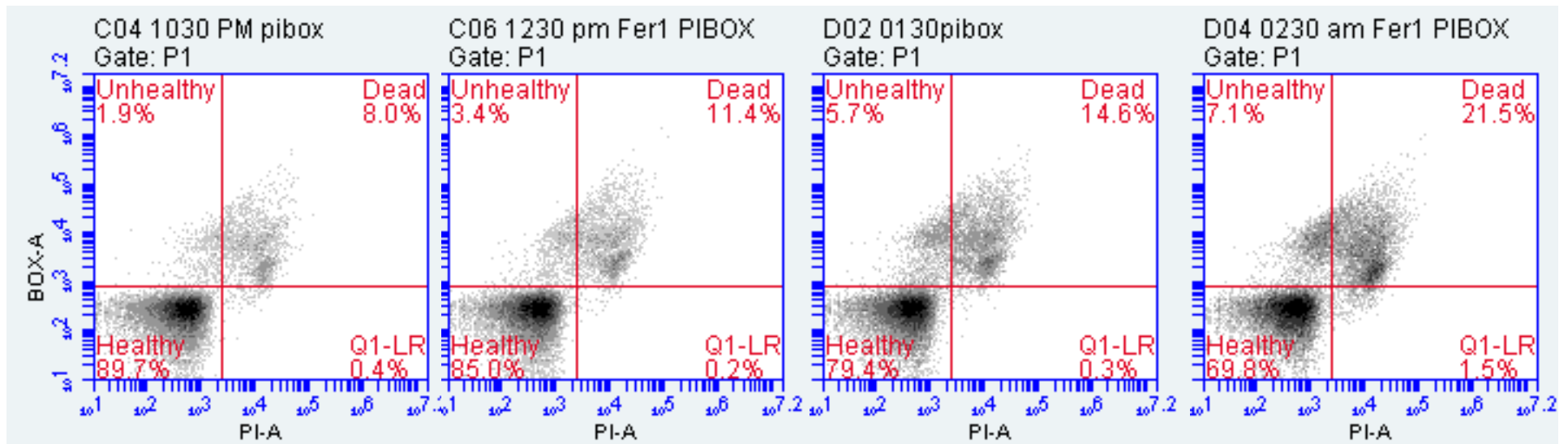


Oxygen consumption & CO₂ evolution



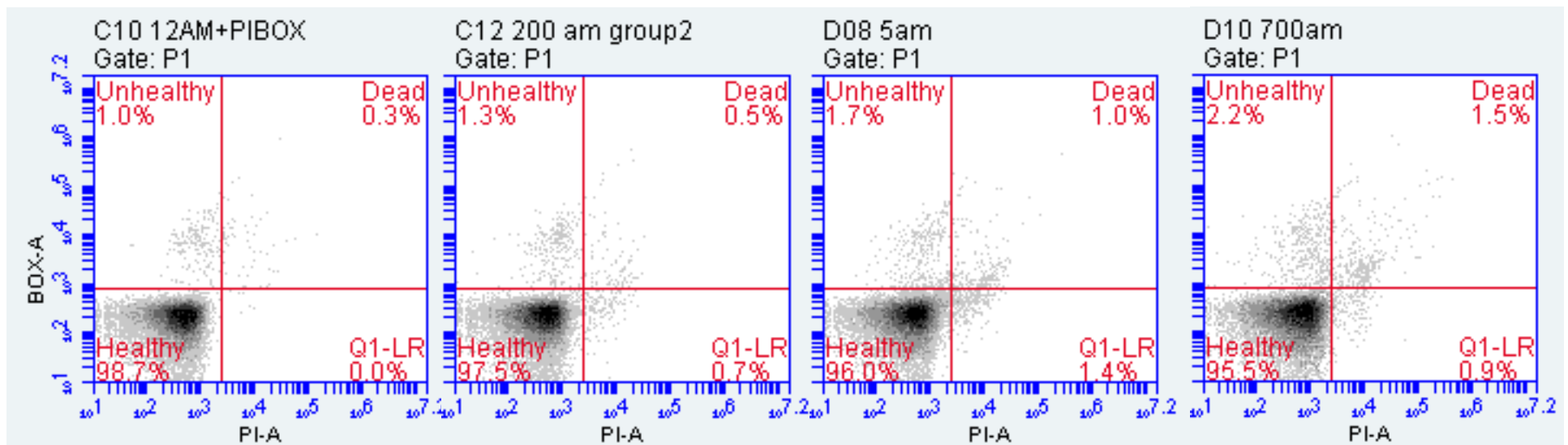
Team 1 – FCM physiology monitoring

*More dead and injured cells –
reflects poor glucose feed control*



Team 2 – FCM physiology monitoring

*More healthy cells – good
glucose feed control*

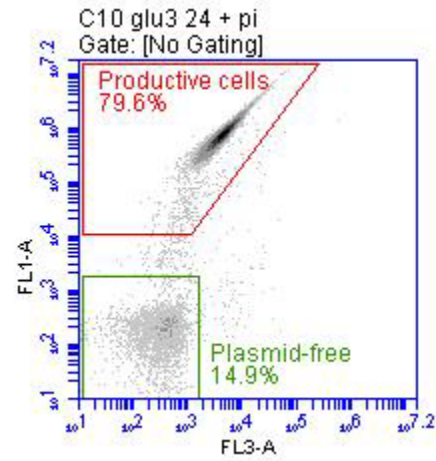
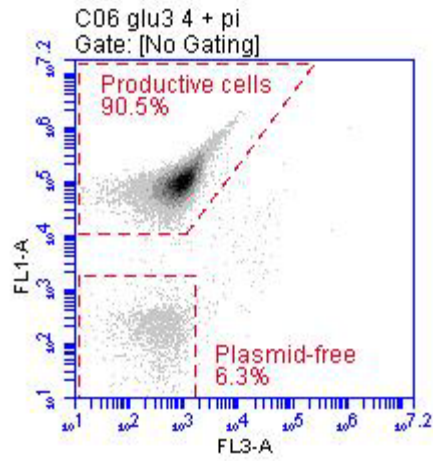
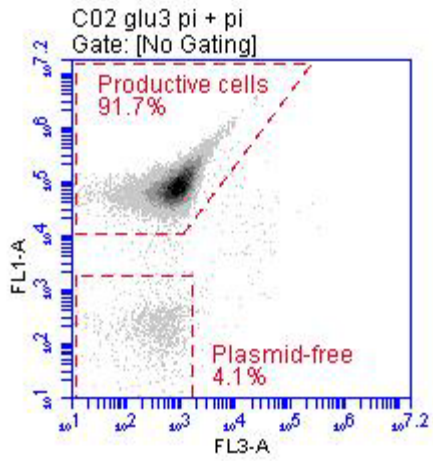


Food microbiology

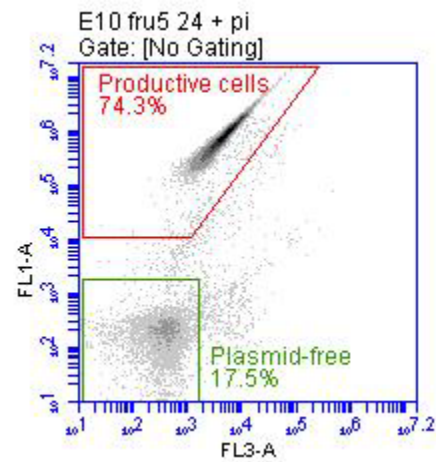
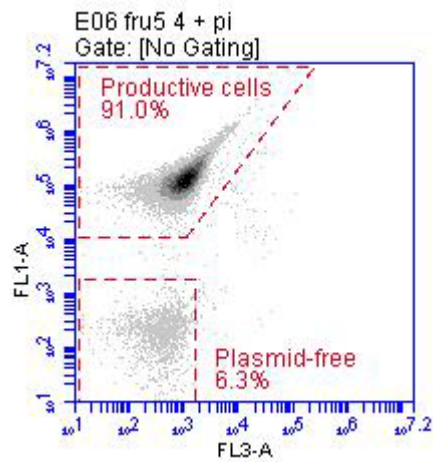
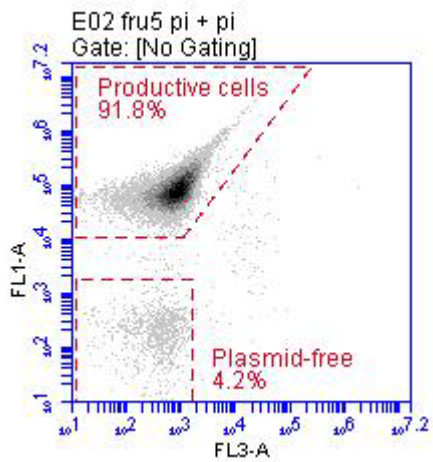
- Use of flow cytometry as a method for monitoring bacteria during food processing
- Acid resistance and acid adaptation
- Differentiation between bacteria and food matrix

MSc / MEng Research projects

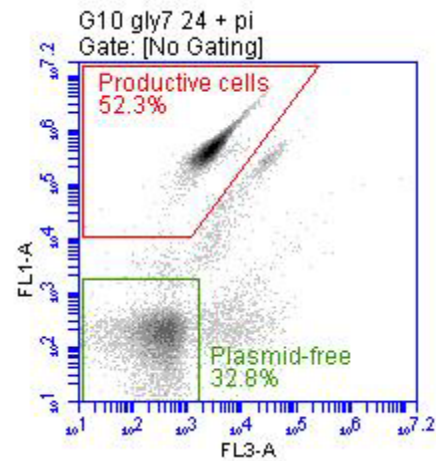
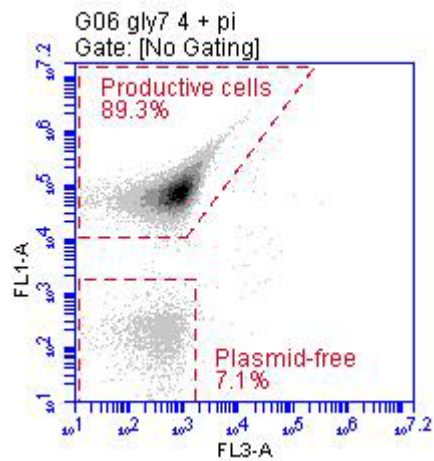
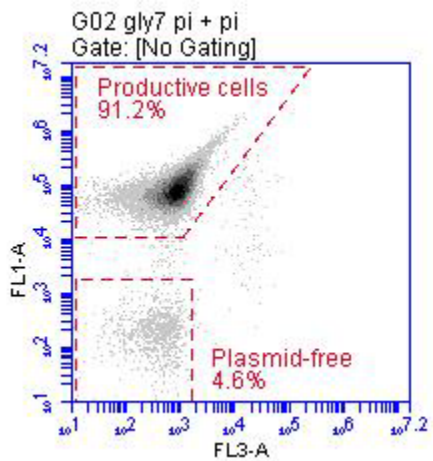
- Monitoring physiology using dyes during biotransformation reactions
- Monitoring poly(3-hydroxybutyrate) (PHB) productivity by bacterial cultures
- Recombinant protein-GFP fusion studies to screen new growth conditions such as choice of carbon source
 - RP-GFP fusions allow measurement of both quantity and folding quality



Glucose



Fructose



Glycerol

Doctoral training

- BBSRC-funded MIBTP (Midlands Integrative Biosciences Training Partnership)
 - Warwick, Birmingham & Leicester Universities
 - 52 PhD students per year
- Extensive training programme:
 - Quantitative skills
 - Placements
 - *Techniques masterclasses*

Doctoral training - master class

- Day 1:
 - Day of lecture / seminar teaching
 - Ability to go into detail and give examples
 - Papers given to students
- Day 2: Reading / preparation
- Day 3:
 - Students give 10 minute talks on papers
 - Practical session

Doctoral training - practical

- *E. coli* wild-type
- *E. coli* GFP⁺
- *S. cerevisiae*
- Comparison of scatter measurements
- Live & Dead cells with a variety of stains (PI, BOX, SYTO9)
- Unknown mixtures (problem-solving aspect)
- Hands-on time is invaluable to apply theory (and test theory)

Challenges and opportunities

- Students have diverse backgrounds
 - Gaps in knowledge (e.g. optics, fluorescence)
 - Students can teach each other
- Students are interested in different areas
 - Comparisons with other techniques
 - Fertile ground for new ideas about methods and measurement techniques
- Students are very keen!

Outlook

- Training in theory of FCM is interdisciplinary and needs life science and physical science knowledge.
- Comparison with previously-used techniques is helpful.
- Combination of taught material with practical experience is very useful.

Acknowledgements

Postdocs: Dr Isaac Vizcaino-Caston & Dr Alfred Fernandez-Castane

PhD students: James Leech, Ikhlaas Kasli, Asma Zulkifly, Hani El Kadri, Hussam Fallatah; Duangkanok Tanangteerapong, Chris Wyre, Amir Anvarian, Louise Hackett

MSc / MEng students: Alifiana Sara, Adna Farah, David Walsh & Christian Mather, Matthew Bridgeman, Mauricio Santos, He Na, Tianqi Wang, Adriana Benarroch, Raúl Mateos González, Ana Álvarez Martín, Shakthiswar Ragu, Kenneth Liu Hung Yaw, Samuel Harvey & David Rothera